

Certificate of Analysis

DIY Human IFN Lambda 3/1/2 (IL-28B/29/28A) ELISA Kit

Catalog No: 61840-1

Lot No: 7867

Expiration: February 17, 2027

| Kit Components | Part No. | Lot No. | Quantity | Amount per vial |
|--------------------|----------|------------|----------|-----------------|
| Capture Antibody | VDY066 | KXU0420081 | 3 | 60 µg |
| Detection Antibody | VDY067 | KLU0920081 | 3 | 24 µg |
| Standard | VDY068 | P395559 | 3 | 80 ng |
| Streptavidin HRP | VDY803 | P330988 | 3 | 2 ml |

Lot Specific Solution Preparations

(Refer to **DIY Human IFN Lambda 3/1/2 (IL-28B/29/28A) ELISA Kit** Protocol for full instructions.)

| Kit Components | Reagents Preparation | Working Concentration |
|--------------------|---|-----------------------|
| Capture Antibody | Reconstitute with 0.5 ml of PBS | 1 µg/ml |
| Detection Antibody | Reconstitute with 1.0 ml of Reagent Diluent | 400 ng/ml |
| Standard | Reconstitute with 0.5 ml of Reagent Diluent | 62.5-4000 pg/ml |
| Streptavidin HRP | Dilute with Reagent Diluent | 40 -fold dilution |

Authorization

Released by: _____ 

Date: April 20, 2026

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.

Rev. 01